

Wortmannin

Phosphatidylinositol 3-kinase Inhibitor

Catalog # tlr1-wtm

For research use only

Version # 15K05-MM

PRODUCT INFORMATION

Content:

- 5 mg Wortmannin

Storage and stability:

- Wortmannin is provided as a translucent film and shipped at room temperature. Store at -20°C. Solid product is stable 1 year at -20°C. Wortmannin is light sensitive. Protect from light.
- Upon resuspension, prepare aliquots of Wortmannin and store at -20°C. Avoid repeated freeze-thaw cycles. Resuspended product is stable for 3 months at -20°C when properly stored.

Quality control:

- Purity: ≥95% (LC)
- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.

DESCRIPTION

Wortmannin is a cell-permeable, fungal metabolite that acts as a potent, selective and irreversible inhibitor of phosphatidylinositol 3-kinase (PI3K)¹. There is increasing evidence of the involvement of PI3K in Toll-like receptor (TLR) signaling². Inhibition of PI3K with wortmannin enhances TLR-mediated inducible nitric-oxide synthase (iNOS) expression, activates NF-κB and up-regulates cytokine mRNA production³. Furthermore, PI3K is required for autophagy⁴. Autophagy is a complex pathway in which cell material can be sequestered and delivered to the lysosome for degradation. Inhibition of PI3K with wortmannin can inhibit autophagic sequestration⁴.

CHEMICAL PROPERTIES

CAS number: 19545-26-7

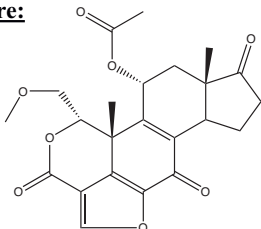
Synonym: KY 12420

Formula: C₂₃H₂₄O₈

Molecular weight: 428.4

Solubility: 10 mg/ml in DMSO

Structure:



METHODS

Preparation of sterile stock solution (20 mM)

1. Add 583 µl DMSO to 5 mg of Wortmannin.
2. Vortex until complete solubilization. Prepare aliquots and store at -20°C. Once Wortmannin is solubilized, dilutions can be prepared using aqueous buffers.

Working concentrations: 100 nM - 10 µM

PROTOCOLS

For reference only; as described in the indicated publications.

Cell Culture Assay³

Cells: Human macrophages

Working concentration: 0.1-1 µM

Incubation time: 15 min

Methods: ELISA, RT-PCR and Western blot

Cell Culture Assay⁴

Cells: Rat hepatocytes

Working concentration: 100 nM

Incubation time: 1 h

Methods: Electron microscopy to study autophagosome formation

Cell Culture Assay⁵

Cells: PANC-1 cells

Working concentration: 200 nM

Incubation time: 1 h

Methods: Electron and fluorescence microscopy, immunoblotting

1. **Arcaro A. and Wymann MP, 1993.** Wortmannin is a potent phosphatidylinositol 3-kinase inhibitor: the role of phosphatidylinositol 3,4,5-trisphosphate in neutrophil responses. *Biochem. J.* 296:297-301. 2. **Fukao T and Koyasu S, 2003.** PI3K and negative regulation of TLR signaling. *Trends Immunol* 24: 358-363. 3. **Hazeki K. et al, 2006.** Opposite effects of wortmannin and 2-(4-Morpholinyl)-8-phenyl-1(4H)-benzopyran-4-one Hydrochloride on Toll-Like Receptor-mediated Nitric Oxide Production: Negative Regulation of Nuclear Factor-κB by Phosphoinositide 3-Kinase. *Mol. Pharmacol.*, 69:1717-4. 4. **Blommaert EF. et al, 1997.** The phosphatidylinositol 3-kinase inhibitors wortmannin and LY294002 inhibit autophagy in isolated rat hepatocytes. *Eur. J. Biochem.* 243: 240-246. 5. **Mancias J. et al, 2014.** Quantitative proteomics identifies NCOA4 as the cargo receptor mediating ferritinophagy. *Nature* 509, 105-9. 6.

RELATED PRODUCTS

Product	Description	Catalog Code
BEZ235	PI3K inhibitor	inh-bez2
LY294002	PI3K inhibitor	tlr1-ly29
Metformin	PI3K inhibitor	tlr1-metf

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Hong Kong: +852 3-622-34-80

E-mail: info@invivogen.com